

Rabbit Anti-Cy7 Antibody, iFluor® 350-Labeled

Catalog Number: 46942

Unit Size: 100 ug

Product Details

Storage Conditions	Refrigerated (2-8 °C), Minimize light exposure
Expiration Date	12 months upon receiving
Concentration	1 mg/mL
Formulation	PBS, BSA

Unit Details

Reconstitution Volume	0.1 mL ddH2O
-----------------------	--------------

Antibody Properties

Clonality	Polyclonal
Host	Rabbit
Immunogen	Cy7 coupled to carrier protein

Biological Properties

Stabilizer	BSA
Appearance	Solid
Preparation	It is produced with Cy7 coupled carrier protein in rabbit and affinity purified with Protein A beads. The antibody is conjugated with iFluor® 350 under optimal conditions.
Soluble In	Water
Application	ELISA, IF, FACS

Spectral Properties

Conjugate	iFluor 350
Excitation Wavelength	345 nm
Emission Wavelength	450 nm

Applications

Rabbit Anti-Cy7 Antibody, iFluor® 350-Labeled is a polyclonal antibody that specifically binds to Cy7-conjugated molecules and is directly labeled with iFluor® 350, a UV-excitable fluorophore with excitation/emission maxima of ~345/450 nm. This direct labeling format supports one-step detection of Cy7-labeled biomolecules without the need for a secondary antibody. The antibody may exhibit cross-reactivity with other near-infrared cyanine dyes. It is recommended for signal amplification, dye conjugation validation, and multiplexed fluorescence assays in microscopy and flow cytometry.